

Anti-Nestin (NES)

Mouse Monoclonal Antibody

Catalog # **N122-561M**

Lot # Z2014-17

Cited Applications

ELISA, IHC

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the Nestin (NES) protein

Cross Reactivity

Human and Rat

Host/Isotype/Clone#

Mouse, IgG

Immunogen

Recombinant human Nestin expressed in bacteria and highly purified. The specific region was amino acids 317 to 630 of the human protein. This is a region which is 55% identical to rat Nestin and 57% identical to mouse Nestin.

Formulation

PBS, pH 7.4 with 0.02% sodium azide.

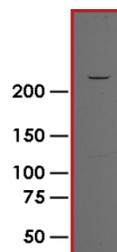
Stability

1yr at -20°C from date of shipment

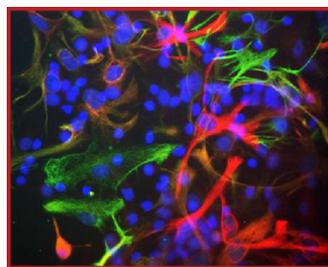
Scientific Background

Nestin (NES) is a Nestin is a Class VI intermediate filament expressed in the developing central nervous system (CNS) in early embryonic neuroepithelial stem cells. This protein has been widely used as a predominant marker for stem / progenitor cells, glioma cells, and tumor endothelial cells in the mammalian CNS. Furthermore, it is a superior angiogenic marker to evaluate neovascularity of endothelial cells in tumor.

Sample Data



Western blot of developing rat brain (P18) homogenate probe with Anti-Nestin (NES) antibody. A single strong band is seen at ~240 kDa.



Mixed cultures of neonatal rat neurons and glia stained with antibody-Nestin (red) chicken antibody to vimentin (green) and DNA (DAPI stain, blue). Astrocytes and neuronal stem cells stain strongly and specifically in a clearly filamentous fashion with the Nestin Antibody. The filamentous staining pattern is as expected as both Nestin and vimentin are components of 10nm filaments. Note that some cells contain Nestin, but do not stain strongly for vimentin and so appear red. Others contain vimentin and not Nestin and so appear green- these are likely to be fibroblastic or endothelial cells. Some cells express both proteins and so appear yellowish. The presence of nestin indicates that the cells are developing astrocytes, neuroblasts or undifferentiated neural stem cells.

Anti-Nestin (NES)

Mouse Monoclonal Antibody

Catalog # **N122-561M**

Lot # **Z2014-17**

Purification **Affinity chromatography**

Stability **1yr at -20°C from date of shipment**

Storage & Shipping **Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.**

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.